



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 2

Catalog Number: C0061-2

Lot Number: A00U0042

Expiration Date: 30 APR 2023

Storage Condition: 2-8°C

Analyte Description	Amount of Analyte in Calibrator	Units	Stock Concentration of Reconstituted Calibrator (Vol = 250 µL)	Units
hu Eotaxin	7 038	pg	28 150	pg/mL
hu Eotaxin-3	23 625	pg	94 500	pg/mL
hu IP-10	11 063	pg	44 250	pg/mL
hu MCP-1	6 850	pg	27 400	pg/mL
hu MCP-4	4 750	pg	19 000	pg/mL
hu MDC	21 875	pg	87 500	pg/mL
hu MIP-1 α	7 013	pg	28 050	pg/mL
hu MIP-1 β	2 300	pg	9 200	pg/mL
hu TARC	3 450	pg	13 800	pg/mL

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

Statement: The above calibrator lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	29 AUG 2018